

Data Sheet

UCHL1

Human, recombinant, N-terminal His-tag

Catalog #: 80351

Lot#: 100826 **Conc.:** 2.38 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
124 mM NaCl, 2.4 mM KCl, 225 mM
imidazole, 10% glycerol, and 3 mM DTT.

Stability: >6 months at -80°C

References:

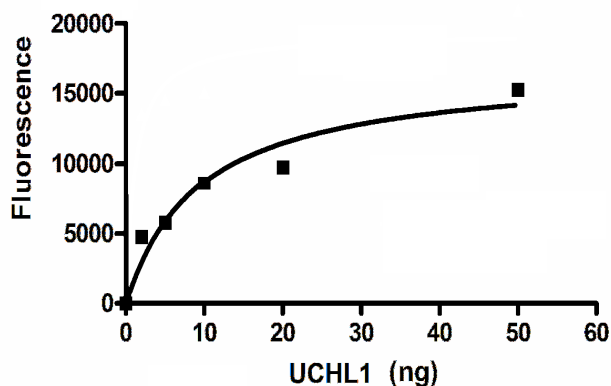
1. Gong, B., *et al.*, (2006). *Cell* **126** (4): 775–788.
2. Healy DG, *et al.*, (2005). *Cell Tissue Res.* **318** (1): 189–94.

Description: UCHL1, also known as Ubiquitin Carboxy-terminal Hydrolase L1 or PGP9.5 (GenBank Accession No. NM_004181), a.a. 2-224(end), with N-terminal His-tag, MW=23.4 kDa, expressed in *E. coli*.

Specific Activity: 7433 pmol/min/μg
Assay Conditions: Reaction was performed in 50 mM Tris, pH 7.4, 1 mM DTT, 0.5 mM EDTA, 500 nM Ub-AMC, and UCHL1. Reaction was incubated at 37°C for 15 min and fluorescent signal was measured at excitation of 340 nm, and emission at 460 nm.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

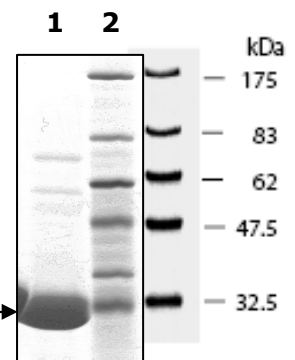


10% SDS-PAGE Coomassie staining

Lane 1:
5 μg UCHL1

Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 23.4 kDa
Purity: >90%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com